

PRODUCT INFORMATION



WAY-208466 (hydrochloride)

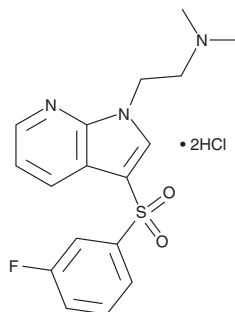
Item No. 11938

CAS Registry No.: 1207064-61-6
Formal Name: 3-[[3-fluorophenyl)sulfonyl]-
N,N-dimethyl-1H-pyrrolo[2,3-b]

MF: C₁₇H₁₈FN₃O₂S • 2HCl
FW: 420.3
Purity: ≥98%

UV/Vis.: λ_{max}: 278 nm
Supplied as: A crystalline solid
Storage: -20°C

Stability: As supplied, 2 years from the QC date provided on the Certificate of Analysis, when stored properly



Laboratory Procedures

WAY-208466 (hydrochloride) is supplied as a crystalline solid. A stock solution may be made by dissolving the WAY-208466 (hydrochloride) in the solvent of choice. WAY-208466 (hydrochloride) is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide, which should be purged with an inert gas. The solubility of WAY-208466 (hydrochloride) in ethanol is approximately 2 mg/ml and the solubility in DMSO and DMF is approximately 25 mg/ml.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. Organic solvent-free aqueous solutions of WAY-208466 (hydrochloride) can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of WAY-208466 (hydrochloride) in PBS, pH 7.2, is approximately 1 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

WAY-208466 is a selective 5-HT₆ receptor agonist (EC₅₀ = 7.3 nM at the human 5-HT₆ receptor).¹ It has been shown to increase cortical GABA levels in rat frontal cortex, producing anxiolytic-like effects *in vivo*.¹

Reference

1. Schechter, L.E., Lin, Q., Smith, D.L., *et al.* Neuropharmacological profile of novel and selective 5-HT₆ receptor agonists: WAY-181187 and WAY-208466. *Neuropsychopharmacology* **33(6)**, 1323-1335 (2008).

WARNING

THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA

This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the complete Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY

Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 11/29/2016

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897
[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM